

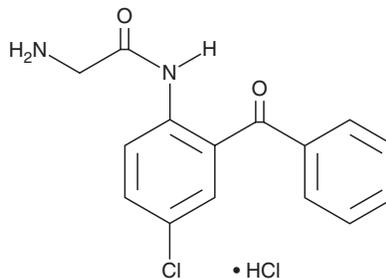
PRODUCT INFORMATION



Nordiazepam Uncyclized Intermediate (hydrochloride)

Item No. 35679

CAS Registry No.: 5504-72-3
Formal Name: 2-amino-N-(2-benzoyl-4-chlorophenyl)-acetamide, monohydrochloride
Synonym: Noravizafone Desglycyl
MF: C₁₅H₁₃ClN₂O₂ • HCl
FW: 325.2
Purity: ≥95%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Description

Nordiazepam uncyclized intermediate (hydrochloride) (Item No. 35679) is an analytical reference standard categorized as a prodrug form of nordiazepam (Item Nos. ISO60174 | 15898).¹ This product is intended for research and forensic applications.

Reference

1. Bundgaard, H. Pro-drugs as drug delivery systems. IX. Reversible cyclization kinetics of 2-aminoacetamido-5-chlorobenzophenone to demethyldiazepam in aqueous solution. *Archiv for Pharmaci og Chemi* **8(1)**, 15-28 (1980).

WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete](#) Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 01/23/2024

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM